
LCK

Product code: 08-170

Full-length human LCK [1-509(end) amino acids of accession number NP_005347.3] was expressed as N-terminal GST-fusion protein (85 kDa) using baculovirus expression system.

Assay platform	: MSA
Substrate	: Srctide
ATP (μ M) Km app / Bin	: 14 / 10
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 1.5
IC50 at 1 mM ATP (nM)	: 14

